

Single-Stranded DNA 7K Ladder Insert

STORE AT -25°C to -15°C UPON RECEIPT

P/N CLS157950: 1 vial of 40 µL of ssDNA 7K Ladder

For Research Use Only, not for use in diagnostic procedures

Properties: Colorless solution, volume per vial: 40 µL, total concentration: 70-110 ng/µL.

Storage buffer components: 50 mM Potassium Acetate, 20mM Tris-acetate, 10mM Magnesium Acetate, 20mM EDTA, ~6% Glycerol, pH ~7.8.

Storage: Store at -25°C to -15 °C. Avoid multiple freeze-thaw cycles. Aliquots of the product may be stored for up to one week at 2°C to 8°C. Do not store in a frost-free freezer.

Handling: Use DNase-free and RNase-free reagents, DNA Low binding tubes, and barrier pipette tips.

Thawing instructions: Thaw just to completion at up to 37°C, vortex a few seconds when fully thawed, and place on ice. To avoid multiple freeze-thaw cycles, make aliquots in DNase-free and RNase-free, DNA Low binding tubes with the required amount typical for daily use. The ssDNA 7K ladder has no obvious stability loss after 3 freeze-thaw cycles.

Shelf life upon receipt: With recommended storage, until expiration specified on the vial.

Ladder gel image on LabChip® GXII Touch™ (P/N CLS138160) using RNA Pico reagent (P/N CLS960012) on the chip of DNA 5K/RNA/CZE HT LabChip® (P/N 760435) using ssDNA 7K.asyx (P/N CLS157952) program.

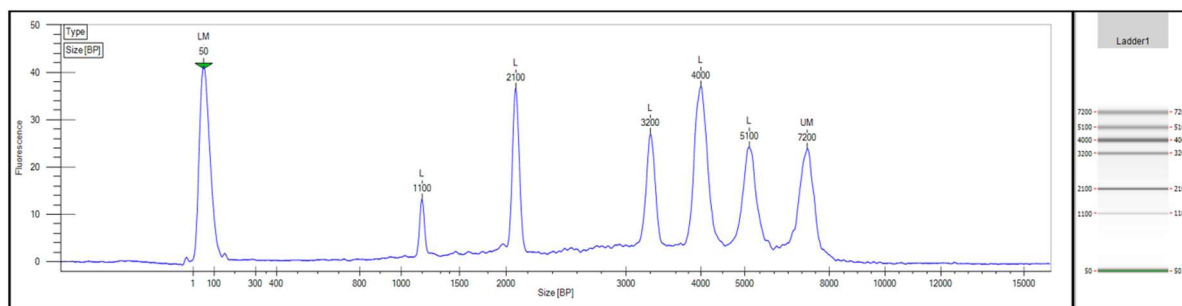


Figure: Example of ladder performance on LabChip® GXII Touch™ with 1:40 dilution by nuclease-free H₂O.